

Rabbit Anti-Human/Mouse/Rat Aven (Apoptosis Caspase Activation Inhibitor)

Instruction Manual

Catalog Number	PK-AB718-2413
Synonyms	Aven Antibody: Apoptosis caspase activation inhibitor
Description	Apoptosis plays a major role in normal organism development, tissue homeostasis, and removal of damaged cells. Disruption of this process has been implicated in a variety of diseases such as cancer. Aven is a recently discovered protein that blocks apoptosis induced by Apaf-1 and caspase-9. It is thought that Aven functions by binding to Bcl-xL, an antiapoptotic member of the Bcl-2 family, and to Apaf-1, possibly interfering with the ability of Apaf-1 to self-associate, suggesting that Aven impedes Apaf-1-mediated caspase activation. Higher levels of Aven mRNA are seen in patients with acute leukemia than in control patients, suggesting that Aven may be useful as a prognostic indicator in leukemia patients. Despite its predicted molecular weight, Aven often migrates at 55 kDa in SDS-PAGE.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Aven antibody was raised against a 14 amino acid peptide from near the carboxy terminus of human Aven.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually. Aven antibody can be used for detection of Aven by Western blot at 1 µg/mL. Despite its predicted molecular weight, Aven often migrates at 55 kDa in SDS-PAGE. Antibody can also be used for immunohistochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Paydas S, Tanriverdi K, Yavuz S, et al. Survivin and aven: two distinct antiapoptotic signals in acute leukemias. <i>Ann. Oncology</i> 2003; 14:1045-50. Chau BN, Cheng EH-Y, Kerr DA, et al. Aven, a novel inhibitor of caspase activation. Binds Bcl-xL and Apaf-1. <i>Mol. Cell</i> 2000; 6:31-40. Lockshin RA, Osborne B, and Zakeri Z. Cell death in the third millennium. <i>Cell Death Differ.</i> 2000; 7:2-7.
Images	Available upon request.
Related Products	Aven Peptide, Cat. No. PK-AB718-2413P Raji Lysate, Cat. No. PK-AB718-1207 Spleen Lysate, Cat. No. PK-AB718-1306

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.